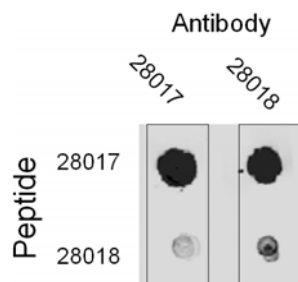


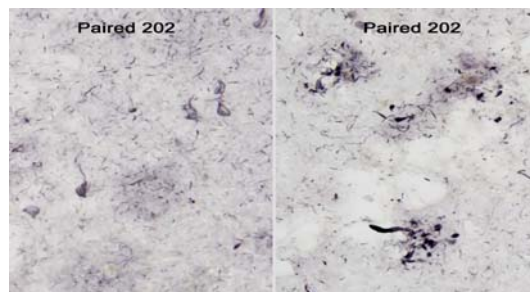


## Product Data Sheet

<b>Product Name:</b>	<b>Anti-Tau (paired202)</b>										
<b>Catalog Number:</b>	28018-025										
<b>Lot Number:</b>	See label on vial										
<b>Product Description:</b>	This polyclonal antibody is supplied as an epitope affinity purified rabbit IgG, 25µg in 125 µl of 1x PBS (pH 7.4) containing 0.05% sodium azide.										
<b>Immunogen:</b>	Rabbit Anti-Tau (paired202) polyclonal antibody was raised against a synthetic peptide (SPGSPGT) corresponding to human Tau around Serine 202.										
<b>Species Reactivity:</b>	Species reactivity includes human, mouse, rat, and bovine, while others remain unknown. This antibody was evaluated for specificity with a dot blot assay (using synthetic Tau peptides) and Western blot assay. By Western blot, immunoreactive bands around 52-68 kDa were detected from mouse brain lysate.										
<b>Application Notes:</b>	The following concentration ranges are recommended starting points for this product. Optimal working concentrations should be determined by the investigator for specific applications.										
	<table> <tr> <td>Dot Blot:</td><td>1:500-2,000</td></tr> <tr> <td>Western Blot:</td><td>1:500-2,000</td></tr> <tr> <td>IHC:</td><td>1:500-2,000</td></tr> <tr> <td>Immunoprecipitation*:</td><td>3.0-5.0 µg/extract from 10<sup>7</sup> cells</td></tr> <tr> <td colspan="2">(*Recommended but not tested)</td></tr> </table>	Dot Blot:	1:500-2,000	Western Blot:	1:500-2,000	IHC:	1:500-2,000	Immunoprecipitation*:	3.0-5.0 µg/extract from 10 <sup>7</sup> cells	(*Recommended but not tested)	
Dot Blot:	1:500-2,000										
Western Blot:	1:500-2,000										
IHC:	1:500-2,000										
Immunoprecipitation*:	3.0-5.0 µg/extract from 10 <sup>7</sup> cells										
(*Recommended but not tested)											



**Figure 1.** Dot blot analysis with phospho- and non-phospho peptides indicated the Anti-Tau (pSer202) antibody (Cat# 28017) to be phospho-peptide specific, while the Anti-Tau (paired202) antibody (Cat# 28018) detected both the phospho- and non-phospho- peptides



**Figure 2.** IHC of the upper layers (II/III) with neurofibrillary tangles (left), and the deeper layers with dystrophic neurites in plaques of entorhinal cortex of Alzheimer disease brains (right), stained with Anti-Tau (paired202) antibody (Cat# 28018), biotinylated secondary antibody avidin-biotin-HRP, visualization by heavy metal intensification of DAB, 40X objective (Courtesy of Dr. Patrick L. McGeer, Kinsmen Laboratory of Neurological Research, University of British Columbia, Vancouver, Canada)



**Figure 3.** Western blot analysis of mouse brain tissue whole cell lysate was probed by Anti-Tau (paired202) antibody (Cat# 28018) at a dilution of 1:500



**Figure 4.** Western blot analysis of non-phosphorylated (1) and phosphorylated (2) recombinant Tau protein probed by Anti-Tau (paired202) antibody (Cat# 28018). Immunoreactive bands were recognized in both non-phospho- and phospho-Tau (Courtesy of Dr. Nichol's Lab, Parkinson's Institute and Clinical Center, Sunnyvale, California)

#### Background:

Tau is a collection of microtubule-associated proteins that is involved in microtubule assembly and stabilization (1). In the adult human brain, six isoforms ranging between 352 and 441 amino acids in length are produced as a result of alternative RNA splicing (2, 3). The expression of Tau isoforms is developmentally regulated, as only the smallest Tau polypeptide is expressed in the fetal brain. Hyper-phosphorylated Tau is the major component of the paired helical filament of Alzheimer's disease. Phosphorylation-dependent anti-Tau antibodies are used to distinguish, between normal brains and Alzheimer's disease brains, specific amino acids that are phosphorylated in Tau. Tau proteins, particularly in developing brains and in Alzheimer brains, were found to phosphorylate *in vivo* at several different sites (4).

#### References:

1. Cleveland, D. et al. *J Mol Biol* **116**, 207 (1977).
2. Goedert, M. et al. *EMBO J.* **8**, 393 (1989).
3. Goedert, M. et al. *Neuron* **3**, 519 (1989).
4. Billingsley, M. et al. *Biochem J* **323**, 577 (1997).

#### Storage:

Store at 2-8°C for up to one year. Avoid repeated freezing and thawing.

#### Related Products:

[Check](#) our website for more Tau antibodies.

This product is for *in vitro* research use only.